

CONFIDENCE IN YOUR DRINKING WATER ANALYSIS WITH GC/MS



Have Confidence in Your Water Testing

About 90% of all drinking water comes from surface and groundwater sources which are susceptible to pollution by the chemicals used in industry and agriculture. Laboratories today are challenged with maintaining compliance with evolving regulations, technologies, and methods. We empower this critical analysis with our GC/MS and Headspace technologies. Coupled with our support expertise, we are a trusted innovator at the forefront of drinking water testing. Have confidence in your water testing with PerkinElmer solutions.

Why the Clarus GC/MS?

When optimizing throughput is key, you can trust us for your most critical volatile organic compound (VOC) requirements and quantification of PAH's, PCB's and pesticides. The Clarus® SQ 8 GC/MS systems outperform regulatory method criteria, and their stability helps you maintain quality control and maximize instrument uptime while ensuring ultra-trace level detection limits.

We stay current with applications and guidance for routine and emerging contaminants in drinking water, delivering solutions for BTEX, fuel oxygenates, and disinfection by-products, as well as for 1,4 dioxane and industrial solvents.



Clarus SQ 8 GC/MS



SMARTsource™

Elite MS Column

What Consumables Do I Need for My Water Analyses?

Maximize the sensitivity of your water analysis by using Elite “MS” GC columns. These phases are designed to be robust for MS applications, with improved thermal stability and ultra-low bleed yielding increased sensitivity.

Facilitate reproducibility in your trace analysis by ensuring an inert flow path for your sample. Use BTO inlet septa and liners from our ultra-inert range, to minimize unwanted sample interactions. Invest in the consumables that are tested and validated to fit your environmental water testing needs.

Select Your Clarus Consumables and Accessories Now:

- ▶ [GC Columns](#)
- ▶ [Injector Septa](#)
- ▶ [Inlet Liners](#)
- ▶ [Syringes](#)
- ▶ [Vials and Caps](#)

www.perkinelmer.com/category/uv-vis-consumables



Your complete solution to get the most out of your analyses, with quality instruments, processes, and data. OneSource instrument services, including comprehensive Instrument Service and Repair, Asset Management and Analytics Services, you have the information you need, in a format you need it, to maximize your lab's productivity.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



Some Water Testing Applications to Suit Your Lab

[The Determination of Low Level Benzene, Toluene, Ethyl Benzene, and Xylenes \(BTEX\) in Drinking Water by Headspace Trap GC/MS](#)

[Determination of 1 ppt of 1,2,3-Trichloropropane and of 1,2-Dibromoethane in Groundwater](#)

[PerkinElmer Partners with Weld County to Optimize Their Water Testing to Include Over 60 Organic Contaminants and a New Method for Methane](#)

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2019, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.